

<b>Product Number:</b>	101361
<b>Lot Number:</b>	503914
<b>Expiration:</b>	28-Mar-2024

Component	Part Number	Lot Number	Control Number	Concentration
SBG Dilution Buffer	100376	316508	316508A	N/A
SBG Concentrate	103397	308712	N/A	50 nM
RGP Reagent	103159	N/A <sup>1</sup>	N/A	N/A

<sup>1</sup> RGP is not Kit Lot Specific

Review/Approval

Sonia Khillan  
Name

QC Manager  
Title

Soni 12/1/23 30 JUN 2023  
Signature/Date